

## Ordering Information

Catalog #	Description
K20-00003	<b>ProteinChip CM Low-Stringency Buffer</b> , 200 ml
C57-30075	<b>ProteinChip CM10 Arrays</b> , A-H format, 12
C50-30011	<b>ProteinChip Cassette-Compatible Bioprocessor</b> , includes ProteinChip array forceps, cassette hold-down frame, 12 blank ProteinChip arrays
C30-00001	<b>ProteinChip CHCA Energy Absorbing Molecules (EAMs)</b> , 5 mg/vial, 20
C30-00002	<b>ProteinChip SPA Energy Absorbing Molecules (EAMs)</b> , 5 mg/vial, 20
C30-00003	<b>ProteinChip EAM-1 Energy Absorbing Molecules (EAMs)</b> , 5 mg/vial, 20

The SELDI process is covered by US patents 5,719,060, 6,225,047, 6,579,719, and 6,818,411 and other issued patents and pending applications in the US and other jurisdictions.

**BIO-RAD**

**Bio-Rad  
Laboratories, Inc.**

*Life Science  
Group*

**Web site** [www.bio-rad.com](http://www.bio-rad.com) **USA** 800 4BIORAD  
**Australia** 61 02 9914 2800 **Austria** 01 877 89 01  
**Belgium** 09 385 55 11 **Brazil** 55 21 3237 9400  
**Canada** 905 712 2771 **China** 86 21 6426 0808  
**Czech Republic** 420 241 430 532  
**Denmark** 44 52 10 00 **Finland** 09 804 22 00  
**France** 01 47 95 69 65 **Germany** 089 318 84 0  
**Greece** 30 210 777 4396  
**Hong Kong** 852 2789 3300  
**Hungary** 36 1 455 8800 **India** 91 124 4029300  
**Israel** 03 963 6050 **Italy** 39 02 216091  
**Japan** 03 5811 6270 **Korea** 82 2 3473 4460  
**Mexico** 52 555 488 7670  
**The Netherlands** 0318 540666  
**New Zealand** 0508 805 500 **Norway** 23 38 41 30  
**Poland** 48 22 331 99 99  
**Portugal** 351 21 472 7700  
**Russia** 7 495 721 14 04 **Singapore** 65 6415 3188  
**South Africa** 27 861 246 723  
**Spain** 34 91 590 5200  
**Sweden** 08 555 12700  
**Switzerland** 061 717 95 55  
**Taiwan** 886 2 2578 7189  
**United Kingdom** 020 8328 2000

# ProteinChip® CM Low-Stringency Buffer

## Instruction Manual

Catalog #K20-00003

For technical support,  
call your local Bio-Rad office, or  
in the US, call **1-800-4BIORAD**  
**(1-800-424-6723)**.

**BIO-RAD**

## Uses

- Protein profiling and biomarker discovery
- Rapid protein analysis to determine purity, mass confirmation, or both

## Introduction

ProteinChip CM low-stringency buffer is designed for use with the ProteinChip CM10 array (catalog #C57-30075). The ProteinChip CM array incorporates a carboxylate chemistry (negatively charged) that acts as a weak cation exchanger. The carboxymethyl (CM) surface binds proteins that are positively charged at a given pH. To control selectivity, the pH of the binding buffer is increased or decreased, depending on the need. By using the low-pH ProteinChip CM low-stringency buffer, an overall net positive charge is imparted on a greater number of proteins within the sample, resulting in more proteins binding. By increasing the pH of the binding and wash buffer, an overall net negative charge is imparted on some of the proteins, resulting in fewer proteins binding, i.e., higher stringency.

## Storage

Store buffer at 2–8°C.

## Buffer Composition

0.1 M sodium acetate, pH 4.0, antimicrobial preservatives, 200 ml — Amount is sufficient to run 12 ProteinChip arrays in a ProteinChip bioprocessor using binding and elution buffer volumes as outlined in the suggested protocol.

## Suggested Protocol

1. Assemble the ProteinChip arrays in the ProteinChip cassette-compatible bioprocessor (catalog #C50-30011).
2. Add 150 µl of ProteinChip CM low-stringency buffer to each well. Vortex for 5 minutes at room temperature.
3. Remove buffer from the wells.
4. Repeat steps 2–3 for a total of two washes.
5. Add 90 µl of ProteinChip CM low-stringency buffer to each well.
6. Add 10 µl of the sample to each well. Vortex for 30 minutes at room temperature.

7. Remove samples from wells.
8. Wash each well with 150 µl ProteinChip CM low-stringency buffer for 5 minutes, with agitation. Repeat twice for a total of three buffer washes.
9. Remove wash buffer from wells and rinse each well with deionized water.
10. Drain wells and remove arrays from the ProteinChip bioprocessor.
11. Allow arrays to air-dry.
12. Apply 1.0 µl ProteinChip energy absorbing molecule (EAM) solution per spot. Two applications of EAM solution can be used in order to increase signal intensity. Allow arrays to air-dry.
13. Analyze the arrays using the ProteinChip SELDI reader.